

## LW-1 Antibody [#C11810]

<b>Catalog Number:</b>	C11810
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9UBD0
<b>Other Names:</b>	Heat shock transcription factor family X-linked 1, LW-1, HSF1, hCG_27154
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from N-terminal of human LW-1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LW-1 antibody detects endogenous levels of total LW-1 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Strausberg R.L., Proc. Natl. Acad. Sci. U.S.A. 99:16899-16903(2002). Wang L., Submitted (APR-1999) to the EMBL/GenBank/DDBJ databases. Mural R.J., Submitted (JUL-2005) to the EMBL/GenBank/DDBJ databases.